

Product Number:	103528
Lot Number:	114116
Expiration:	21-May-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26×10^9 Beads/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 15 Jul 2021

Signature/Date